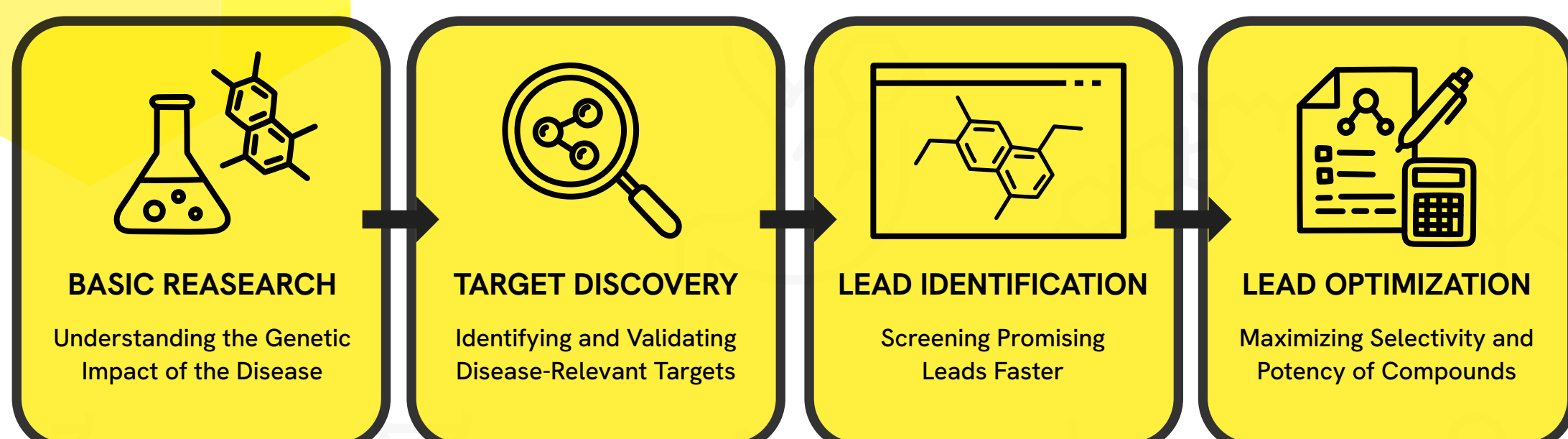
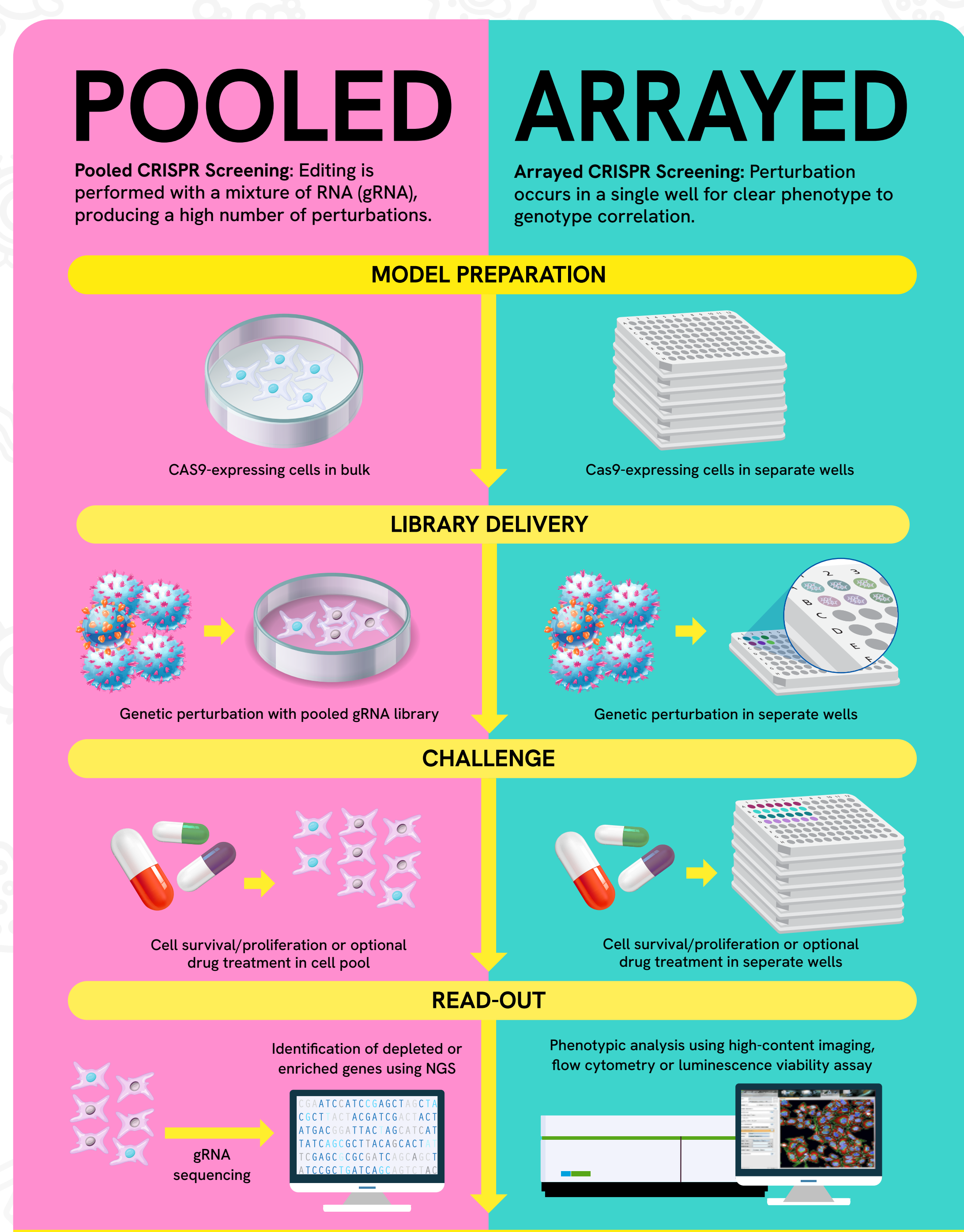


# Functional genomic screening in drug discovery

Functional genomic screens enable modulation of hundreds or thousands of genes in a single experiment to identify genetic pathways, cellular processes, novel therapeutic targets, and to genetically profile existing or potential therapeutics.

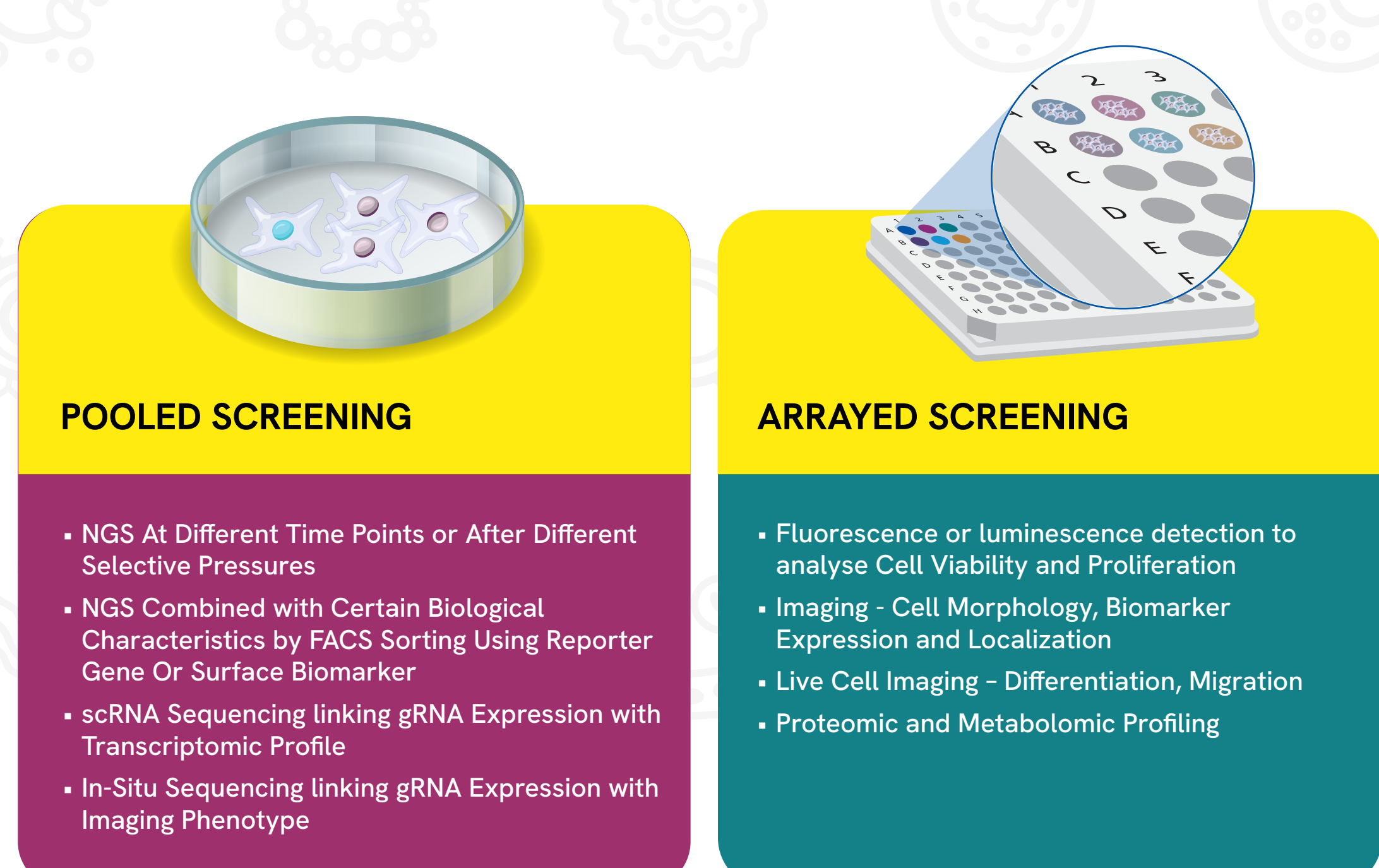


Gene expression can be regulated using either RNAi or CRISPR technology, where RNAi represses gene expression at the mRNA level (knockdown), while CRISPR works at the DNA level and can permanently knockout or modulate genes. The two CRISPR screening approaches are:

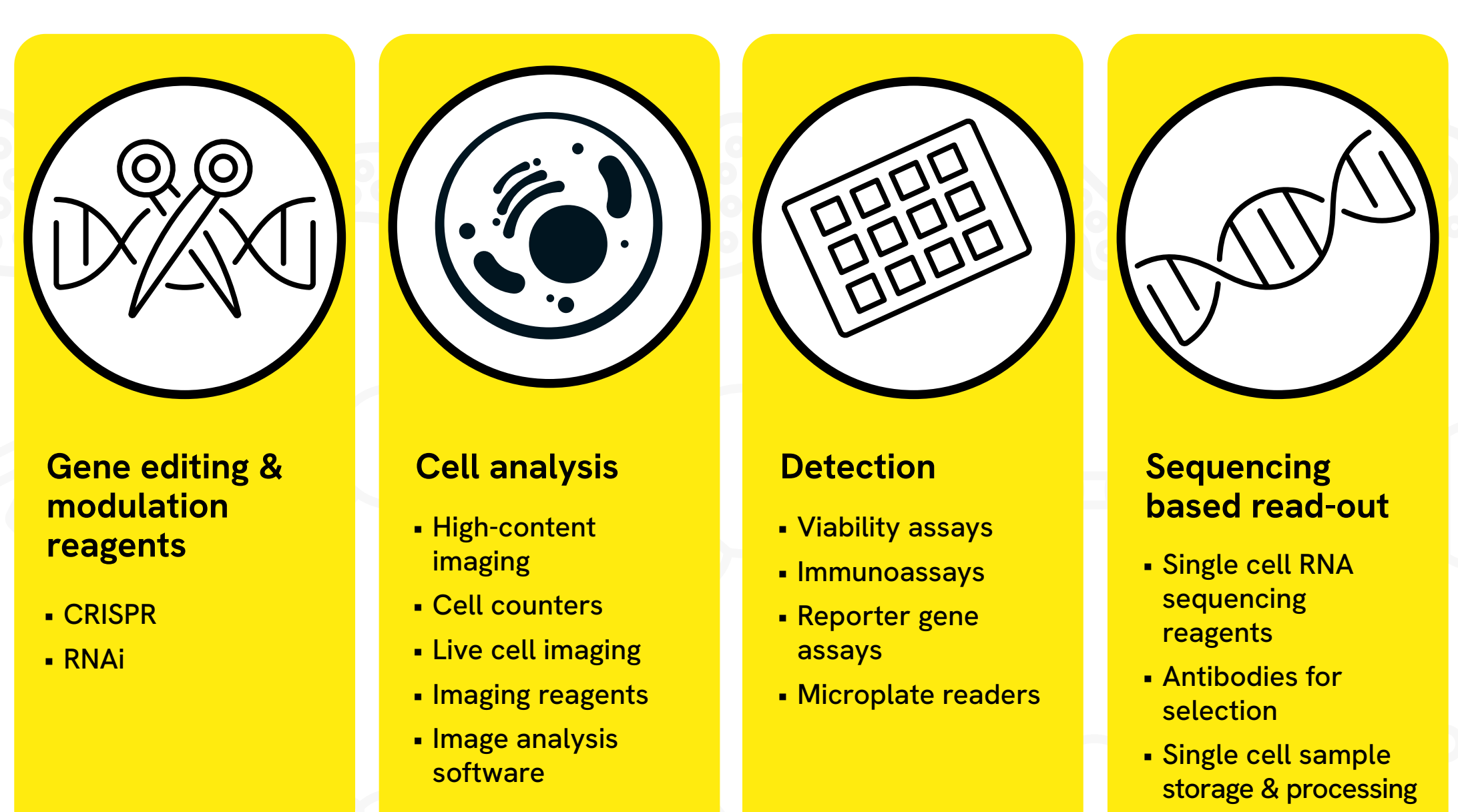


|               | POOLED SCREENING  | ARRAYED SCREENING   |
|---------------|---|---|
| Advantages    | <ul style="list-style-type: none"> <li>Less expensive</li> <li>Large number of genes can be integrated at once</li> <li>Suitable for longer <i>in vivo</i> studies</li> <li>No specialized/automated equipment needed</li> </ul>  | <ul style="list-style-type: none"> <li>Multi-parametric read-outs</li> <li>Suitable for more complex cell models</li> <li>More choices in reagent formats, including non-viral transfection and cherry-picking options</li> <li>Clear phenotype to genotype correlation</li> <li>Improved editing efficiency by targeting the same gene with several gRNAs in one well</li> </ul> |
| Disadvantages | <ul style="list-style-type: none"> <li>Only simple read-outs: cell death or survival, reporter gene and sortable biomarkers</li> <li>An edited cell type can influence other cells in the mixed population negatively</li> <li>Very large cell population needed at start and to be maintained in large flasks</li> </ul> | <ul style="list-style-type: none"> <li>Lab automation equipment needed</li> <li>Assays often suffer from low transfection/electroporation efficiencies</li> </ul>   |

## Read-out methods



## Products



## Functional genomic screening services

Comprehensive solutions to perform pooled and arrayed screens for you with our CRISPR knockout, CRISPR activation, CRISPR inhibition, and dual CRISPRa/i screening platforms.

[LEARN MORE](https://www.revivity.com/)

<https://www.revivity.com/>

revvity